

MHC-II Monoclonal Mouse Antibody (IPO-10)

Product Description

MAB IPO-10 defines the antigen that appears on B cell progenitors following HLA-DR and preceding CD10, CD19, CD22, CD37 and cym. It is expressed on resting B cells and then reappears and persists in cytoplasm and on cell surface until cytoplasmic Ig appears. It is a useful antibody for diagnostics of neoplasms of B cell origins. It reacts with human B cell lines Daudi, Raji, Namalva, EB-3, RPMI-8226 (50% of cells). The MAb does not label T cell lines, blood granulocytes, thymocytes or bone marrow stromal fibroblasts. No significant changes are detected after PHA or ConA stimulation while LPS and PWM stimulated cultures after 18-48h show decreased number of antigen-positive cells but in final terms of cultivation antigen is expressed again. This MAb labels B cell leukemias and some lymphomas. Hairy cell leukemia strongly reacts and 70% of B cell CLL and some B-NHL were also positive. IPO-10 reacts with AMML cells and in a majority of Hodgkin's disease cases a significant percentage of affected lymph node cells were detected.

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

Stock status: Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 1090, Anti-MHC-II (IPO-10)

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

1. Anticancer Res (1988) 8(6):1377-1380.
2. Tissue Antigens (1989) 33: 151.
3. Neoplasma (1991) 39(1):3-9.

Product attributes

Antibody number	#1090
Antibody reactivity (target)	MHC-II
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	IPO-10
Isotype	IgG3, kappa
Molecular weight	36 kDa
Synonyms	Ia3; Ia-3; H-2Ea; H2-Ea; A1323765; I-Ealpha; E-alpha-f; MHC-H2-Ea
Human gene symbol	HLA-DRA
Entrez gene ID	3122
SwissProt	P01903
Unigene	520048
Immunogen	Spleen cells of a patient with hairy cell leukemia (Daudi cells)
Antibody target cellular	Plasma membrane
Species reactivity	Human, Monkey
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	Ramos, Daudi or HuT78 cells. Tonsil or lymph node
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Antibody research areas	Immunology, MHC antigens

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and of LI-COR Bioscience.

This datasheet was generated on August 30, 2024 at 06:19:38 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-mhc-ii-hla-dra-ipo-10/>